

Certificate of Registration

QUALITY MANAGEMENT SYSTEM - ISO 9001:2008

This is to certify that:

Bio-Rad Laboratories Limited
Bio-Rad House
Maxted Road
Hemel Hempstead
HP2 7DX
United Kingdom

Holds Certificate Number:

FM 635742

and operates a Quality Management System which complies with the requirements of ISO 9001:2008 for the following scope:

Sales, distribution and after-sales service of life science research products, medical diagnostic products and special spectral databases and chemistry publishing software.

For and on behalf of BSI:



Frank Lee, EMEA Compliance & Risk Director

Original Registration Date: 27/01/1994

Latest Revision Date: 26/08/2015

Effective Date: 26/08/2015

Expiry Date: 25/08/2018



Page: 1 of 3

...making excellence a habit.™

Certificate No: FM 635742

Location	Registered Activities
Bio-Rad Laboratories Limited Bio-Rad House Maxted Road Hemel Hempstead HP2 7DX United Kingdom	Sales, distribution and after-sales service of life science research products, medical diagnostic products and special spectral databases and chemistry publishing software.
DiaMed (GB) Limited Unit 2 Whitehill Business Centre Dalkeith Midlothian EH22 2QB United Kingdom	Sales, distribution and after-sales service of life science research products, medical diagnostic products and special spectral databases and chemistry publishing software.
Bio-Rad Laboratories BV Winninglaan 3 Temse 9140 Belgium	Sales, distribution and after-sales service of life science research products, medical diagnostic products and special spectral databases and chemistry publishing software.
Bio-Rad Laboratories AB Solna Strandväg 3 17154 Solna Sweden	Sales, distribution and after-sales service of life science research products, medical diagnostic products and special spectral databases and chemistry publishing software.
Bio-Rad Laboratories Torgbygget Nydalsveien 33 N-0421 Oslo Norway	Sales, distribution and after-sales service of life science research products, medical diagnostic products and special spectral databases and chemistry publishing software.
Bio-Rad Laboratories OY Valimotie 1A FIN-00380 Helsinki Finland	Sales, distribution and after-sales service of life science research products, medical diagnostic products and special spectral databases and chemistry publishing software.
Bio-Rad Laboratories Symbion Science Park Fruebjergvej 3 2100 Kobenhavn O Denmark	Sales, distribution and after-sales service of life science research products, medical diagnostic products and special spectral databases and chemistry publishing software.
DiaMed Benelux, Winninglaan 3 9140 Temse Belgium	Sales, distribution and after-sales service of life science research products, medical diagnostic products and special spectral databases and chemistry publishing software.

Original Registration Date: 27/01/1994

Effective Date: 26/08/2015

Latest Revision Date: 26/08/2015

Expiry Date: 25/08/2018

Page: 2 of 3

This certificate was issued electronically and remains the property of BSI and is bound by the conditions of contract.

An electronic certificate can be authenticated [online](#).

Printed copies can be validated at www.bsigroup.com/ClientDirectory

Certificate No: FM 635742

Location	Registered Activities
Bio-Rad Laboratories SA-NV Winninglaan 3 9140 Temse Belgium	Sales, distribution and after-sales service of life science research products, medical diagnostic products and special spectral databases and chemistry publishing software.
DiaMed Fennica OY Valmotie 1A FIN-00380 Helsinki Finland	Sales, distribution and after-sales service of life science research products, medical diagnostic products and special spectral databases and chemistry publishing software.



Original Registration Date: 27/01/1994

Latest Revision Date: 26/08/2015

Effective Date: 26/08/2015

Expiry Date: 25/08/2018

Page: 3 of 3

This certificate was issued electronically and remains the property of BSI and is bound by the conditions of contract.
An electronic certificate can be authenticated [online](#).
Printed copies can be validated at www.bsigroup.com/ClientDirectory

Information and Contact: BSI, Kitemark Court, Davy Avenue, Knowlhill, Milton Keynes MK5 8PP. Tel: + 44 845 080 9000
BSI Assurance UK Limited, registered in England under number 7805321 at 389 Chiswick High Road, London W4 4AL, UK.
A Member of the BSI Group of Companies.